



# MA100H Moisture Analyzer

Fully Automatic Infrared Moisture Analyzer with Halogen Heater

Moisture Analyzer



- Time-saving and convenient alternative to a drying oven
- Comprehensive range of options for flexible adaptation to various drying tasks
- Ideally suited for ambitious measuring tasks in the fields of quality control or research
- Utmost safety due to self-calibrating system and full GMP suitability
- Ideally suited for samples with high moisture content due to quick control behaviour

Max. weighing capacity	100 g
Repeatability, typical	Starting at an initial sample weight of approx. 1 g: $\pm 0.1\%$   starting at an initial sample weight of approx. 5 g: $\pm 0.02\%$
Readability	1 mg, 0.01%   0.1 mg, 0.001%
Typical sample quantity	5 – 15 g
Reading display	% moisture content (optionally available with conversion factor)   % dry weight   RATIO value   mg weight loss   g residue   g/kg residue   g/l residue
Temperature range and settings	30°C–180°C, stand-by temperature selectable from 30–100°C in increments of 1 degree
Sample heating	Infrared heating using a halogen round heater
Heating programs	Standard, quick, gentle and phase drying
Shutoff criteria	Optional: <ul style="list-style-type: none"> <li>- Fully automatic</li> <li>- SPRM mode (optimization to a reference system)</li> <li>- Semi-automatic (1–50 mg/5–300 sec.)</li> <li>- Semi-automatic (0.1–5.0%/5–300 sec.)</li> <li>- Timer setting (3x 0.1–999 min.)</li> <li>- Timer setting + fully/semi-automatic (2x 0.1–999 min. + automatic)</li> <li>- Manual</li> </ul>
Access to sample chamber	Motorized cover
Measuring program	30 programs saved in a non-volatile memory (freely selectable parameters)
Data storage	Result statistics for the last 9999 measurements/program
Operator guidance features	Menu-driven, alphanumeric dialogue text (5 languages selectable)
Parameter   data entry	<ul style="list-style-type: none"> <li>- Optimized text entry via softkeys</li> <li>- Numeric entry via keypad</li> <li>- Optional data entry via barcode scanner</li> </ul>
Report printout	<ul style="list-style-type: none"> <li>- Printout using the optional, internal printer YDP01MA</li> <li>- Configurable GLP-report for measuring results   calibration   adjustment (5 languages selectable)</li> </ul>



Easy to clean



With optional printer YDP01MA

Test equipment monitoring	<ul style="list-style-type: none"> <li>- Internal calibration weight</li> <li>- Automatic reproTEST for the weighing system</li> <li>- Automatic temperature calibration with optional temperature adjustment set YTM03MA</li> <li>- Documentation in accordance with the applicable quality assurance requirements</li> </ul>
Safety	Parameter settings password-protected against unauthorized access
Voltage frequency	48-60 Hz
Power consumption	Max. 700 VA
Operating temperature range	10°C...30°C
Housing dimensions (W × D × H)	350 × 453 × 156
Weight	Approx. 8.0 kg

#### Available Models

	Supply voltage
MA100H-000115V1	115 V +15%, -20%
MA100H-000230V1	230 V +15%, -20%

#### Accessories

6965542	Disposable sample pans, 80 units, aluminum, round, Ø 90 mm
6906940	Fiberglass pad, Ø 90 mm for pasty and fatty samples, hard quality, 80 units
6906941	Fiberglass pad, Ø 90 mm for liquid and fatty samples, soft quality, with high suction force, 200 units
YAT01MA	Single-use pipettes for weighing liquid samples, 500 units
YSC02	SartoCollect, software for data communication between moisture analyzer and PC (connecting cable must be ordered separately) Prerequisite: MS Windows 2000   Windows XP Professional, Vista Ultimate, Windows 7 Ultimate.
YCW4528-02	External calibration weight, 50 g (E2) with DKD certificate
YTM03MA	Temperature adjustment set with manufacturer certificate for adjustment of the heat source
YDP01MA	Data printer, suitable for integration
YDP20-OCE	Data printer, suitable for external verification, with date, time, statistics, numerator including cable
6906918	Ink ribbon for data printer
6906937	Paper rolls for data printer, 5 rolls, 50 m each

#### Equipment Supplied

- Moisture analyzer with data interface
- Power cord
- User manual
- 80 alu sample pans
- In-use dust cover for keypad
- 1 pair of forceps
- Laminated short-form manual (6 languages)

Sartorius Weighing Technology GmbH  
Weender Landstrasse 94-108  
37075 Goettingen, Germany

Phone +49.551.308.0  
Fax +49.551.308.3289

info.mechatronics@sartorius.com  
www.sartorius-mechatronics.com

Subject to change without notice.  
Printed in Germany on paper that has been bleached  
without any use of chlorine. | W  
Publication No.: W--2030-e11081  
Order No.: 98649-012-20